



EpiCypher 2023  
CONFERENCE

[EPICYPHER.COM/EPI2023](https://epicypher.com/epi2023)

# EPICYPHER 2023 SPONSORSHIP OPPORTUNITIES

NOVEMBER 5-10TH, 2023  
CANCÚN, MEXICO



# SPONSORSHIP OPPORTUNITIES

## RECOGNITION LEVELS & BENEFITS

### BRONZE

- Your company logo linked to your website on EpiCypher 2023 Conference Webpage.
  - Your company logo linked to your website included in the EpiCypher 2023 Digital Program.
  - Your company logo on sponsor slide which runs between each session.
  - Individual slide advertisement to appear during slide loop.
- 

### SILVER

- Your company logo linked to your website on EpiCypher 2023 Conference Webpage.
  - Your company logo linked to your website included in the EpiCypher 2023 Digital Program.
  - Your company logo on sponsor slide which runs between each session.
  - Individual slide advertisement to appear during slide loop.
  - Attendee registration summary, which lists all registrants and their contact information.
  - All-inclusive attendee registration for one person.
  - Showcase your company with a table-top exhibit, at the entrance to meeting hall where coffee breaks will be held.
- 

### GOLD

- Your company logo linked to your website on EpiCypher 2023 Conference Webpage.
  - Your company logo linked to your website included in the EpiCypher 2023 Digital Program.
  - Your company logo on sponsor slide which runs between each session.
  - Individual slide advertisement to appear during slide loop.
  - Attendee registration summary, which lists all registrants and their contact information.
  - All-inclusive attendee registration for one person.
  - Showcase your company with a table-top exhibit, at the entrance to meeting hall where coffee breaks will be held.
  - Gift bag inserts distributed to all attendees.
  - Your company logo will be placed on select EpiCypher2023 Conference emails and social media posts.
- 

### CO-SPONSOR (1 AVAILABLE)

- Your company logo linked to your website on EpiCypher 2023 Conference Webpage.
- Your company logo linked to your website included in the EpiCypher 2023 Digital Program.
- Your company logo on sponsor slide which runs between each session.
- Individual slide advertisement to appear during slide loop.
- Showcase your company with a table-top exhibit, at the entrance to meeting hall where coffee breaks will be held.
- Attendee registration summary, which lists all registrants and their contact information.
- Gift bag inserts distributed to all attendees.
- Your company logo will be placed on select EpiCypher2023 Conference emails and social media posts.
- All-inclusive attendee registration for three people.
- Co-Sponsor logo placement on EpiCypher 2023 Conference Webpage and Program.

# SPONSORSHIP BREAKDOWN

BENEFITS	BRONZE	SILVER	GOLD	CO-SPONSOR
Your company logo linked to your website on EpiCypher 2023 Conference Webpage.	X	X	X	X
Your company logo linked to your website included in the EpiCypher 2023 Digital Program.	X	X	X	X
Your company logo on sponsor slide which runs between each session.	X	X	X	X
Table-top exhibit		X	X	X
Individual slide advertisement to appear during slide loop.		X	X	X
Attendee List		X	X	X
Gift bag insert			X	X
Your company logo will be placed on select EpiCypher2023 Conference emails and social media posts.			X	X
Co-Sponsor logo placement on all EpiCypher 2023 Conference materials.				X
All-inclusive attendee registration		X	X	X
All-inclusive attendee registration				X
All-inclusive attendee registration				X
<b>Cost</b>	<b>\$2,500</b>	<b>\$7,500</b>	<b>\$10,000</b>	<b>\$20,000</b>

## À LA CARTE ITEMS

Table-Top Exhibit	\$3,000
1 Attendee All-Inclusive Registration	\$2,400
Poster Session Sponsor (2 Available)	\$5,000

Interested in taking the next step?  
 Contact us at  
[epicypher2023@epicypher.com](mailto:epicypher2023@epicypher.com)